



8-(6-Aminoethyl)-amino-ATP

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-807S	50 µl (10 mM)
NU-807L	5 x 50 µl (10 mM)

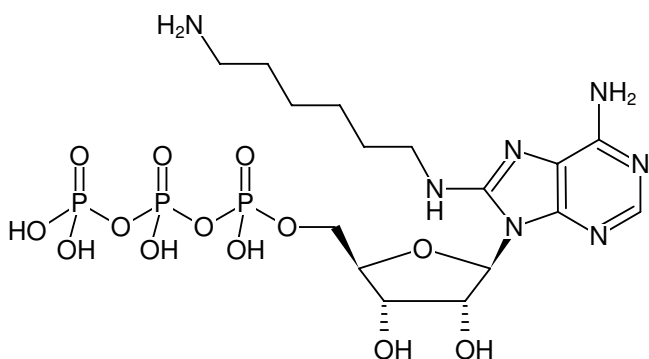
Specific Ligands:

Agonist for P2Y receptor^[1]

Selected References:

[1] Brunstock *et al.* (1994) Structure activity relationships for derivatives of adenosine-5'-triphosphate as agonists at P2 purinoreceptors: heterogeneity within P2X and P2Y subtypes. *Drug Dev. Res.* **31** (3):206.

Komori *et al.* (2008) Measurement system for simultaneous observation of myosin V chemical and mechanical events. *BioSystems* **93**:48.



Structural formula of 8-(6-Aminoethyl)-amino-ATP

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₆H₃₀N₇O₁₃P₃ (free acid)

Molecular Weight: 621.37 g/mol (free acid)

Exact Mass: 621.11 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 273 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)